

PRODUCT INFORMATION



Thromboxane B₂ MaxSpec[®] Standard Item No. 10007237

CAS Registry No.: 54397-85-2

Formal Name: 9 α ,11,15S-trihydroxythromba-5Z,13E-dien-1-oic acid

Synonym: TXB₂

MF: C₂₀H₃₄O₆

FW: 370.5

Purity: \geq 95%

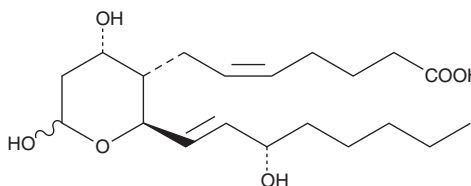
Supplied as: A solution in methyl acetate; in a deactivated glass ampule

Concentration: 100 μ g/ml (nominal); see certificate of analysis for verified concentration

Storage: -20°C

Stability: \geq 3 years; *Stability testing is ongoing to ensure concentration accuracy. The certificate of analysis and product expiry date will be updated upon completion of testing.*

Special Conditions: Store upright and unopened at -20°C. Warm to room temperature prior to opening. Light sensitive.



Description

Thromboxane B₂ (TXB₂) is a non-enzymatically derived, stable, inactive metabolite of TXA₂, which is highly unstable.¹⁻³ Serum levels of TXB₂ positively correlate with platelet COX-1 activation.^{4,5} Urinary levels of TXB₂ reflect intrarenal TXA₂ synthesis, while its metabolites, 11-dehydro TXB₂ (Item No. 19500) and 2,3-dinor TXB₂ (Item No. 19510), reflect systemic TXA₂ secretion.^{2,6,7}

TXB₂ MaxSpec[®] standard is a quantitative grade standard of TXB₂ (Item No. 19030) that has been prepared specifically for mass spectrometry or any application where quantitative reproducibility is required. The solution has been prepared gravimetrically and is supplied in a deactivated glass ampule sealed under argon. The concentration was verified by comparison to an independently prepared calibration standard. This TXB₂ MaxSpec[®] standard is guaranteed to meet identity, purity, stability, and concentration specifications and is provided with a batch-specific certificate of analysis. Ongoing stability testing is performed to ensure the concentration remains accurate throughout the shelf life of the product.

Note: The amount of solution added to the vial is in excess of the listed amount. Therefore, it is necessary to accurately measure volumes for preparation of calibration standards. Follow recommended storage and handling conditions to maintain product quality.

References

1. Needleman, P., Moncada, S., Bunting, S., *et al.* Identification of an enzyme in platelet microsomes which generates thromboxane A₂ from prostaglandin endoperoxides. *Nature* **261**(5561), 558-560 (1976).
2. Patrono, C., Ciabattini, G., Pugliese, F., *et al.* Estimated rate of thromboxane secretion into the circulation of normal humans. *J. Clin. Invest.* **77**(2), 590-594 (1986).
3. Uyama, O., Matsumoto, M., Fujisawa, A., *et al.* Plasma concentrations of thromboxane B₂ in patients with hypertension or cerebrovascular disease. *Prostaglandins Med.* **7**(3), 199-207 (1981).
4. Arantes, F.B.B., Menezes, F.R., Franci, A., *et al.* Influence of direct thrombin inhibitor and low molecular weight heparin on platelet function in patients with coronary artery disease: A prospective interventional trial. *Adv. Ther.* **37**(1), 420-430 (2020).
5. Ferroni, P., Riondino, S., Vazzana, N., *et al.* Biomarkers of platelet activation in acute coronary syndromes. *Thromb. Haemost.* **108**(6), 1109-1123 (2012).
6. Lawson, J.A., Patrono, C., Ciabattini, G., *et al.* Long-lived enzymatic metabolites of thromboxane B₂ in the human circulation. *Anal. Biochem.* **155**(1), 198-205 (1986).
7. Patrono, C., Ciabattini, G., Patrignani, P., *et al.* Evidence for a renal origin of urinary thromboxane B₂ in health and disease. *Adv. Prostaglandin Thromboxane Leukot. Res.* **11**, 493-498 (1983).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/01/2021

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM